

R-Phycoerythrin Cyanine 5.0

Product Number	D14-000
Lot Number	XXXX
Concentration:	10.47 mg/mL
Amount	2mg
Store at 2-8°C	-

<u>Form/Shipping & Storage</u> Supplied as liquid. Upon receipt, store at 2-8° C. DO NOT FREEZE.

<u>Handling</u>

We recommend that the investigator determine the appropriate working concentration for their specific application. Light Sensitive!

<u>Buffer</u>

100mM Sodium Phosphate (pH 7.0) + 5mM EDTA and 0.5% sodium azide

<u>Stability</u> Product should be stored at 2-8°C in the dark.

Note

For research use only, not for diagnostic or therapeutic use.

General Information on APC:

• Approximate molecular weight250,000 kDa• Subunit structure $(\alpha\beta)_6^{\Box}$ • Absorbance/excitation maximum565>540 >498 >650• Emission maximum664 nm• 664/570 Emission ratio4.95

4985 Winchester Blvd. Frederick, MD 21703 301.732.5415 www.columbiabiosciences.com For technical inquiries: info@columbiabiosciences.com For sales inquiries: sales@columbiabiosciences.com